

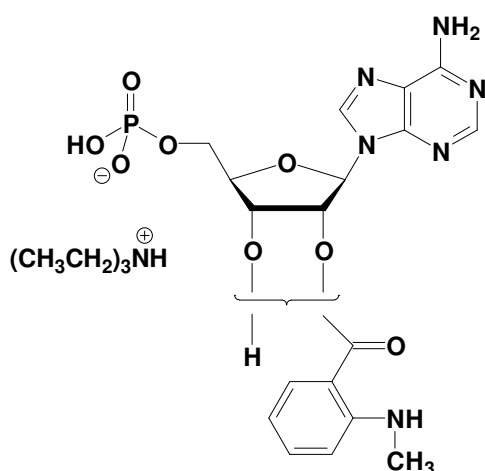
## mant-AMP

### 2'/3'-(N-Methyl-anthraniloyl)-adenosine-5'-monophosphate, Triethylammonium salt

Cat. No.	Amount
NU-236S	150 Units
NU-236L	750 Units

#### Selected References:

Hiratsuka T. *et al.* (1983) New Ribose-modified Fluorescent Analogs of Adenine and Guanine Nucleotides Available as Substrates for Various Enzymes. *Biochim Biophys Acta* **742** :496.



**Cat. No.:** NU-236

**Molecular Formula:** C<sub>18</sub>H<sub>21</sub>N<sub>6</sub>O<sub>8</sub>P (Anion)

**Molecular Weight:** 479.40 (Anion)

**Purity:** > 95%, clear aqueous solution, pH 7.5

#### Spectroscopic properties:

$\lambda_{\max}$  255/355 nm;  $\epsilon$  23300/5800;  $\lambda_{\text{exc}}$  355nm;  
 $\lambda_{\text{em}}$  448 nm

#### Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

#### For research use only!

1 unit = 1  $\mu$ l of a 10 mM solution